



11526 Sorrento Valley Rd SteA2
San Diego Ca 92121
Tel: 1.858.202.1401
1.858.829.3082
Fax: 1.858.259.5734
Email: info@bpsbioscience.com

Data Sheet

PDE2A

human, recombinant, N-terminal GST tag
Catalog #: 60020

Formulated in: 25 mM Tris-HCl, pH 8.0,
100 mM NaCl, 0.05% Tween-20, 50% glycerol,
and 3 mM DTT.

Stability: >6 months at -80°C

References:

1. Seybold, J., et al., Blood 105 (9), 3569-3576 (2005).
2. Martinez, S.E., et al., Proc. Natl. Acad. Sci. U.S.A. 99 (20), 13260-13265 (2002).

Description:

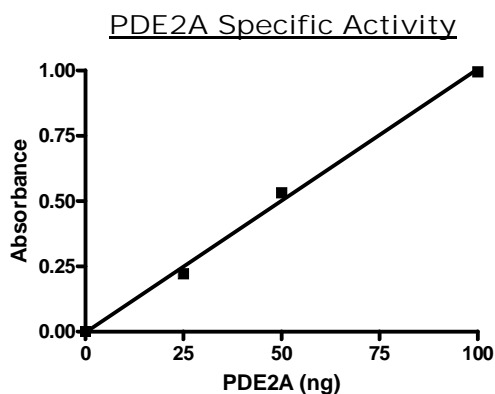
Human PDE2A1 (GenBank Accession No. NM_002599), full length, with N-terminal GST tag, MW=132 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: 2080 pmol/min/μg. 1 unit is defined as the amount of enzyme that will convert 1 pmole of 3', 5'-cAMP to 5' AMP per min at 30 °C in a reaction buffer. Assay conditions: 10 mM Tris-HCl, pH7.4, 10 mM MgCl₂, 1 mM MnCl₂, 200 μM cAMP, 2.5 kU 5' nucleotidase, 37°C, 20 min.

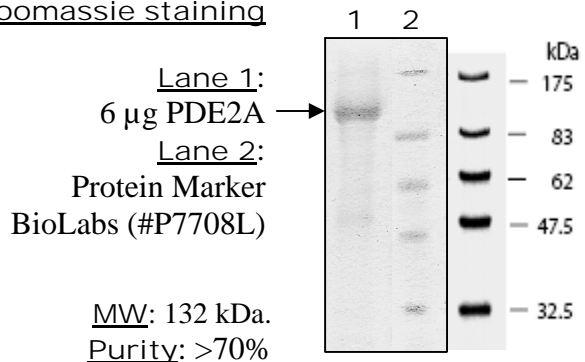
Application:

Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance



10% SDS-PAGE
Coomassie staining



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** or **1.858.829.3082**, Fax **1.858.259.5734**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com